

TREVIGEN[®]

MATERIAL SAFETY DATA SHEET

PRODUCT NAME: **HT Chemiluminescent PARG Assay Kit**

CATALOG NUMBER: **4682-096-K**

Hazards identified with this product are those associated with the following component(s). Refer to the material safety data sheet for the listed item(s).

<u>Catalog Number</u>	<u>Name</u>
4668-050-01	PARP Enzyme
4671-096-02	20X PARP Buffer
4671-096-03	10X PARP Cocktail
4680-096-01	PARG Enzyme
4680-096-02	10X PARG Buffer
4671-096-06	10X Activated DNA
4678-096-P	Histone-coated strip wells
4680-096-03	6,9-diamino-2-ethoxyacridine in DMSO
4680-096-04	Strep-HRP80
4675-096-01	PeroxyGlow [™] A
4675-096-02	PeroxyGlow [™] B

Warning Label

Irritant.

Irritating to eyes, respiratory system and skin.

In case of contact with eyes, rinse immediately with plenty of water and seek medical advice.

Wear suitable protective clothing, gloves, and eye/face protection.

The physical, chemical, and toxicological properties of the product(s) may not yet have been fully characterized. Trevigen recommends the use of gloves, lab coats and eye protection while using or handling any reagent(s). Trevigen assumes no liability for damage resulting from handling or contact with the above product(s).